

**TNP-UTP**

2',3'-O-Trinitrophenyl-uridine-5'-triphosphate, Triethylammonium salt

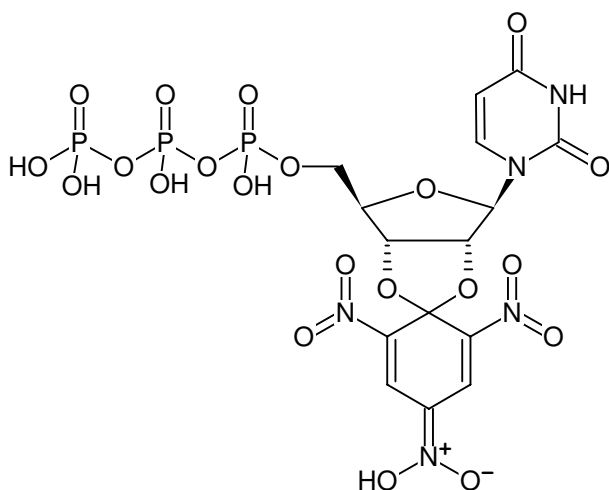
Cat. No.	Amount
NU-226S	100 µl (10 mM)
NU-226L	5 x 100 µl (10 mM)

**Selected References:**

Suryanarayana *et al.* (2009) Differential Inhibition of Various Adenylyl Cyclase Isoforms and Soluble Guanylyl Cyclase by 2',3'-O- (2,4,6-Trinitrophenyl)-Substituted Nucleoside 5'-Triphosphates. *J. Pharmacol. Exp. Ther.* **330** (3):687.

Goettle *et al.* (2007) Molecular analysis of the interaction of Bordetella pertussis adenylyl cyclase with fluorescent nucleotides. *Molecular Pharmacology* **72** (3):526.

Sprang *et al.* (2006) Broad Specificity of Mammalian Adenylyl Cyclase for Interaction with 2',3'-Substituted Purine- and Pyrimidine Nucleotide Inhibitors. *Mol. Pharmacol.* **70**:878.



Structural formula of TNP-UTP

**For research use only!****Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>15</sub>H<sub>16</sub>N<sub>5</sub>O<sub>21</sub>P<sub>3</sub>**Molecular Weight:** 695.23 g/mol (free acid)**Exact Mass:** 694.96 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** orange**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 262/408/470 nm,  
 ε 10.2/26.4/18.5 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5), λ<sub>exc</sub> 408/470 nm,  
 λ<sub>em</sub> 552 nm